

Rhodamine Phalloidin Reagent ab235138

23 References

概述

产品名称	Rhodamine Phalloidin试剂
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Cell-based (qualitative)
产品概述	Rhodamine Phalloidin Reagent ab235138 is one of a series of phalloidin conjugates that are used to stain actin filaments in fluorescence microscopy. The tetramethyl rhodamine dye can be easily detected with a fluorescent microscope at Ex/Em = 546/575 nm.

Phalloidin conjugates are convenient probes for labeling, identifying and quantitating animal or plant actin filaments in formaldehyde-fixed and permeabilized tissue sections, cell cultures or cell-free experiments.

Review other popular phalloidin dye conjugates, including [Phalloidin-iFluor 488](#), [Phalloidin-iFluor 647](#), [Phalloidin-iFluor 594](#), and [Phalloidin-iFluor 555](#), search the website to see [all phalloidin conjugates](#), or read the [phalloidin staining protocol](#).

说明	<p>Staining fixed cell or tissue samples with phalloidin conjugates is very simple; it requires a single 20-90 min incubation with the phalloidin, followed by 3 short wash steps. Phalloidin staining can be combined with antibody-based staining by adding the phalloidin conjugate during either the primary or secondary antibody incubation step.</p> <p>When used in unfixed samples, phalloidin binding leads to a decrease in the disassociation rate of actin subunits from the ends of actin filaments, essentially stabilizing actin filaments through the prevention of filament depolymerisation.</p>
----	---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

平台	Fluorescence microscope
----	-------------------------

性能

存放说明	Store at -20°C. Please refer to protocols.
------	--------------------------------------------

组件	300 tests
Rhodamine Phalloidin Conjugate	1 x 300 tests

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors